

GLOBAL PROFICIENCY TESTING PROGRAM FOR HUMAN BONE MARROW COLONY ASSAYS Session-Specific Information

Session	October 2011
Sample ID	BM27

For a complete step-by-step video that demonstrates how to receive the kit components and set up the CFC assay for the Proficiency Testing Program, please visit our website at www.stemcell.com/proficiencyvideo.

STORAGE AND HANDLING

- Frozen bone marrow samples should be stored at **-135°C or colder** (i.e. ultra-low freezer or liquid nitrogen). If cells are not stored appropriately, the viability and CFC potential of the cells will be compromised.
- MethoCult™ GF and IMDM +2% FBS should be stored at **-20°C**.
- Remaining kit components can be stored at room temperature (15 - 25°C).

PREPARATION OF CELL DILUTION

10X Cell Dilution Concentration	2 x 10 ⁵ cells per mL
--	----------------------------------

- The 10X cell dilution concentration is specific to each cell sample. It is required for calculating the dilution of the thawed sample with IMDM + 2% FBS prior to plating, as described on the Proficiency Testing Program Data Submission Worksheet (Catalog #29168) that was included in this shipment.

DEADLINE FOR INITIATING CFC ASSAY

Thaw and plate cells by	November 14, 2011
--------------------------------	-------------------

DEADLINE AND INSTRUCTIONS FOR SUBMITTING RESULTS

Submit results by	November 29, 2011
--------------------------	-------------------

Submit results in one of three ways:

- **Online** with the Proficiency Testing Data Submission Forms available on our website at www.proficiencytesting.com
- **E-Mailing** the completed Proficiency Testing Program Data Submission Worksheet (Catalog #29196) to proficiency@stemcell.com
- **Faxing** the completed Proficiency Testing Program Data Submission Worksheet (Catalog #29196) to 1.604.877.0704 or 1.800.567.2899 (North America) (Attention Education Department)

WARNING

This kit contains human bone marrow cells. The bone marrow was obtained from a normal donor who was screened for Hepatitis B surface (HBs) antigen, Hepatitis Be antigen, Hepatitis Be antibody, Hepatitis C antibody and HIV-1 antibody. However, the sample cannot be guaranteed free of pathogenic agents. Human cells should be considered potentially infectious and handled accordingly.

Copyright © 2011 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design and MethoCult are trademarks of STEMCELL Technologies Inc.

28455 Lot #11A38324
October 2011